

**Purified Mouse Anti-human/mouse
AHNAK1 Antibody *EM-09, monoclonal***

Catalog number: V103060

Unit size: 0.1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, mouse
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	AHNAK1
Clone	EM-09

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	WB, IHC(F), ICC, IP

Applications

Neuroblast differentiation-associated protein AHNAK is a protein with a molecular weight of 735 kDa, located in the actin cytoskeleton, costamere and nucleus of cells. Sequencing of neuroblast differentiation-associated protein AHNAK has demonstrated it contains a primary structural unit, the PDZ domain. Neuroblast differentiation-associated protein AHNAK reacts with cadherin, RNA and identical protein. It is an integral part of voltage-gated calcium channel activity and RNA splicing.